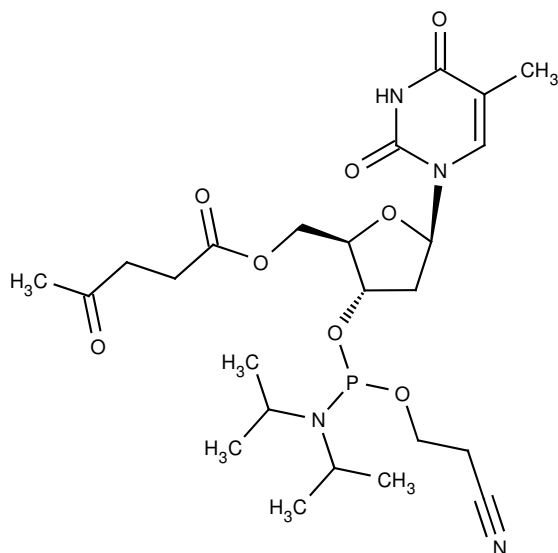


5'-Levulinyl-3'-CEP Thymidine

Cat. No.	Amount
N-ANP-5004-250	250 mg
N-ANP-5004-500	500 mg
N-ANP-5004-1G	1 g
N-ANP-5004-2G	2 g



Molecular Formula: C₂₄H₃₇N₄O₈P

Molecular Weight: 540.55 g/mol

Storage conditions: store at -20°C

Shelf life: 12 months

Purity: >98% (HPLC)

Appearance: thick oily form

For research use only!